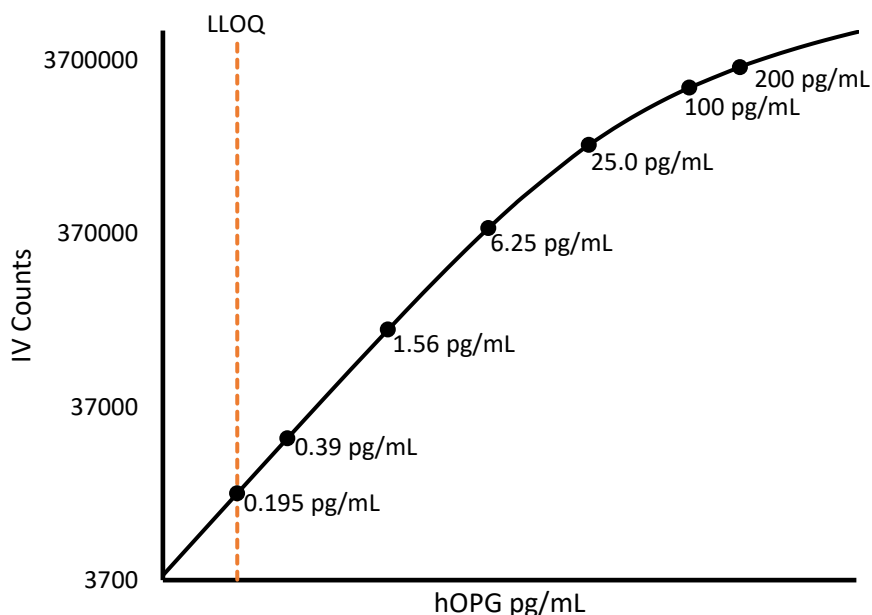


Description – OPG (Osteoprotegerin)

OPG (Osteoprotegerin), also known as OCIF (osteoclastogenesis inhibitory factor), is a 60 kDa protein that plays an important role in bone density and homeostasis by inhibiting osteoclasts and promoting osteoclast apoptosis. OPG binds to RANKL, preventing the ligand from causing bone resorption. Levels of OPG in blood have been associated with several diseases including atherosclerosis, cardiovascular disorders, juvenile Paget’s, and osteoporosis.

Calibration Curve: Calibrator concentrations and Lower Limit of Quantification are depicted in the figure below. This standard curve is for demonstration purposes; end users should prepare a standard curve for each assay run.



Minimum Required Dilution (MRD)

Diluted Sample volume (1:50 Dilution)*	50 µL per measurement
---	-----------------------

*See Kit Instructions for details

Endogenous Serum and Plasma Readings: Healthy EDTA plasma samples (n=3) were measured.

% Above LOD	100%
% Above LLOQ	100%

Assay Range: The upper end of the dynamic range is equal to the top calibrator concentration multiplied by MRD.

Analytical LLOQ	0.195 pg/mL
Functional LLOQ (x MRD)	9.75 pg/mL
LOD	17.5 fg/mL
Assay Range	0 – 10,000 pg/mL

Note: Data described were developed during assay development. Under different assay conditions, assay may perform differently than shown. For complex matrices such as serum or plasma, assay diluent optimization (for example by adding blocking agents) may improve performance of these matrices in this assay.